

HTRF® Research Reagents

HTRF cAMP Gi kit - 1K pts

Product Number : 62AM9PEB Lot Number : 06Q

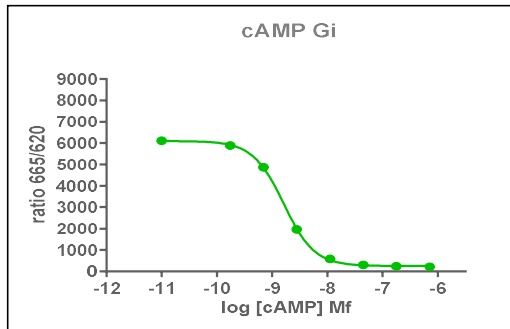
Expiration date : 2024-11-02 Document version : 1

Reagents :	Lot n°
HTRF Lysis & Detect. Buf. 8 - 13 mL	06RA
HTRF Stimulation Buf. 1 (5X) - 8 mL	24RA
HTRF cAMP Gs Dynamic & Gi kit - Std	17AR
HTRF cAMP Gi kit - 1K pts Eu Cryptate reagent	06RA
HTRF cAMP Gi kit - 1K pts d2 antibody	06RA
/	/
/	/

Results :

Incubation : 1h at RT
 Plate : 384 wells small volume white plate 20 µL
 Reader : Pherastar Plus
 Fitting model : nonlinear regression, sigmoidal dose response (variable slope)

cAMP	Conc. (nM)	DF %
Std 0	0	2784
Std 1	0,69	2683
Std 2	2,78	2200
Std 3	11,12	829
Std 4	44,5	174
Std 5	178	44
Std 6	712	12
Std 7	2848	1



Assay characteristics :

	Found value	Specifications
S/B	29	27-40
Hillslope	-1,556	-1.35-2.04
IC50 Nm	1,604	1.25-3
blanc k 620 mm	30156	

HTRF cAMP Gi kit - 1K pts - Lot n° : 06Q

Accepted

Rejected

Quality control date : 2023-12-27	Supervisor's initials : MP
-----------------------------------	----------------------------

The information provided in this document is valid for the specified lot number and date of analysis. This information is for reference purposes only and does not constitute a warranty or guarantee of the product's suitability for any specific use. Revvity, Inc., its subsidiaries, and/or affiliates (collectively, "Revvity") do not assume any liability for any errors or damages arising from the use of this document or the product described herein. REVVITY EXPRESSLY DISCLAIMS ALL WARRANTIES, INCLUDING WARRANTIES OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE, REGARDLESS OF WHETHER ORAL OR WRITTEN, EXPRESS OR IMPLIED, ALLEGEDLY ARISING FROM ANY USAGE OF ANY TRADE OR ANY COURSE OF DEALING, IN CONNECTION WITH THE USE OF INFORMATION CONTAINED HEREIN OR THE PRODUCT ITSELF.